



# Solvents

## Application Note

Environmental

### Authors

Agilent Technologies, Inc.

### Introduction

Fast gas chromatography with an Agilent PoraPLOT U column separates ten solvents in five minutes.



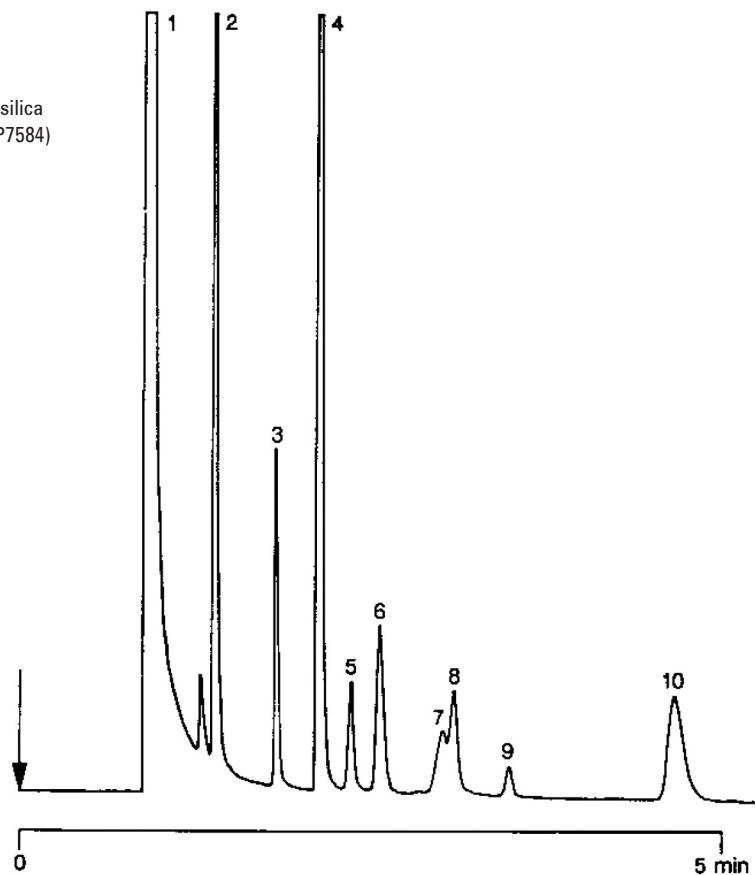
**Agilent Technologies**

## Conditions

Technique : GC-Wide-bore  
Column : Agilent PoraPLOT U, 0.53 mm x 25 m fused silica  
PLOT PoraPLOT U (df = 20 µm) (Part no. CP7584)  
Temperature : 150 °C  
Carrier Gas : H<sub>2</sub>, 80 kPa (0.8 bar, 11.4 psi), 12 mL/min  
Injector : Direct  
Detector : TCD  
Sample Size : 100 - 1000 ppm

## Peak identification

1. air
2. methanol
3. ethanol
4. acetone
5. isopropanol
6. hexane
7. tert. butylalcohol
8. propanol
9. ethyl acetate
10. heptane



[www.agilent.com/chem](http://www.agilent.com/chem)

This information is subject to change without notice.

© Agilent Technologies, Inc. 2011

Printed in the USA

31 October, 2011

First published prior to 11 May, 2010

A00681



**Agilent Technologies**